



CERTIFICATE OF ANALYSIS

Page 1 of 1

Product Description: Calibrator 1

Catalog Number: C0060-2

Lot Number: A00U0272

Expiration Date: 31 MAY 2027

Storage Temperature: 2-8°C

Intended Assay Species: Human

Concentration Assignment:

Analyte	Amount of Analyte in Calibrator	Unit of Measure	Stock Concentration of Reconstituted Calibrator	Unit of Measure
IFN- γ	38375	pg	153500	pg/mL
IL-1 β	6038	pg	24150	pg/mL
IL-2	2650	pg	10600	pg/mL
IL-4	2488	pg	9950	pg/mL
IL-6	2450	pg	9800	pg/mL
IL-8	2900	pg	11600	pg/mL
IL-10	5075	pg	20300	pg/mL
IL-12p70	7300	pg	29200	pg/mL
IL-13	2238	pg	8950	pg/mL
TNF- α	3763	pg	15050	pg/mL
IL-5	5650	pg	22600	pg/mL
GM-CSF	13625	pg	54500	pg/mL
IL-17A	38875	pg	155500	pg/mL
VEGF-A	4038	pg	16150	pg/mL

Reconstitution Volume: 250 μ L

To calculate the **Calibrator Standard 1 (top of curve) concentration**, please refer to the dilution-factor used in the **Prepare Calibrator Standards** section of the respective product insert. Divide the stock concentration of the reconstituted calibrator by this dilution factor to generate the working calibrator standard 1 concentration.

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		08 AUG 2022